

HTRF® Research Reagents

# HTRF Human IFN alpha kit - 500 pts

Product Number : 62HIFNAPEG Lot Number : 06H

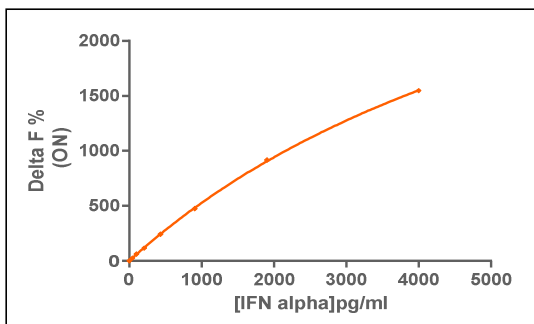
Expiration date : 2024-12-21 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	19G
HTRF Detect. Buf. 3 - 1.5 mL	28A
HTRF Human IFN alpha kit - Std	05RA
HTRF Human IFN alpha kit - 500 pts d2 antibody	05A
HTRF Human IFN alpha kit - 500 pts Eu Cryptate antibody	06RA

**Results :**

Incubation : **2h at RT**  
 Plate : **384 wells small volume white plate 20 µL**  
 Reader : **Pherastar Plus**  
 Fitting model : **4 Parameter Logistic (4PL 1/y2)\* model**

hIFN alpha	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	47	29%	22-32
Std 2	98	55%	43-66
Std 3	206	110%	106-122
Std 4	432	221%	199-277
Std 5	907	449%	386-537
Std 6	1905	892%	765-961
Std 7	4000	1530%	1247-1551



**Assay characteristics :**

	Found value	Specifications Rev 01
Detection limit	7.2 pg/ml	< 47 pg/ml

HTRF Human IFN alpha kit - 500 pts - Lot n° : 06H

Accepted

Rejected

Quality control date : 2023-11-29	Supervisor's initials : PL
-----------------------------------	----------------------------

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.